

Mitochondria Isolation Kit for Tissue ab110168

★★★★★ [2 Abreviews](#) [58 References](#) [1 图像](#)

概述

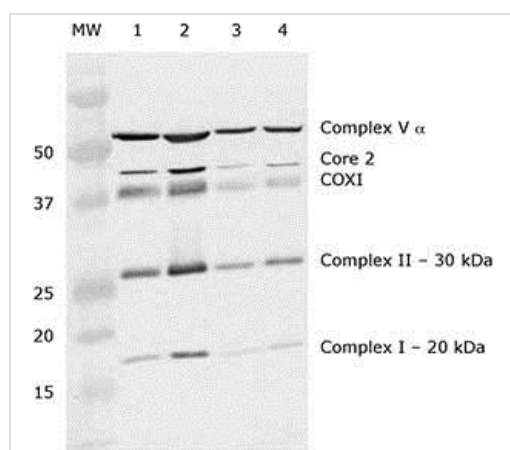
产品名称	Mitochondria分离试剂盒for组织
样品类型	Tissue
检测时间	1h 00m
产品概述	<p>ab110168 (MS850) contains sufficient reagents for up to 10 benchtop mitochondria isolations from hard or soft tissues. Isolated mitochondria can be analyzed by Western Blotting, Immunocapture, or Activity assay.</p> <p>Mitochondrial isolation protocol summary:</p> <ul style="list-style-type: none">- suspend minced tissue in isolation buffer- homogenize using dounce homogenizer- spin at 1,000 g for 10 min and retain supernatant- spin at 12,000 g for 15 min and retain pellet- resuspend pellet in isolation buffer, spin at 12,000g for 15 min, collect pellet and repeat spin- resuspend pellet in isolation buffer and analyze or freeze
说明	Tested tissues: brain, liver and heart

性能

存放说明 Store at +4°C. Please refer to protocols.

组件	1 kit
Isolation Buffer	1 x 100ml
Wash Buffer	1 x 30ml

图片



Western blot - Mitochondria Isolation Kit for Tissue
(ab110168)

Rat liver mitochondria were isolated using Abcam's **ab110169/MS851** Mitochondria Isolation Kit for Tissue (with Dounce Homogenizer). 2 µg (lane 1) and 10 µg (lane 2) of mitochondria were tested with Abcam's **ab110413/MS604** (MitoProfile® Total OXPHOS Rodent WB Antibody Cocktail). Mitochondria are enriched and so Western blot band intensity of mitochondrial proteins is greater in these mitochondrial preparations when compared to homogenate at respective amounts, see lane 3 and 4.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors